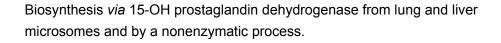
15-Ketoeicosa-6E,8Z,11Z,13Etetraenoic acid



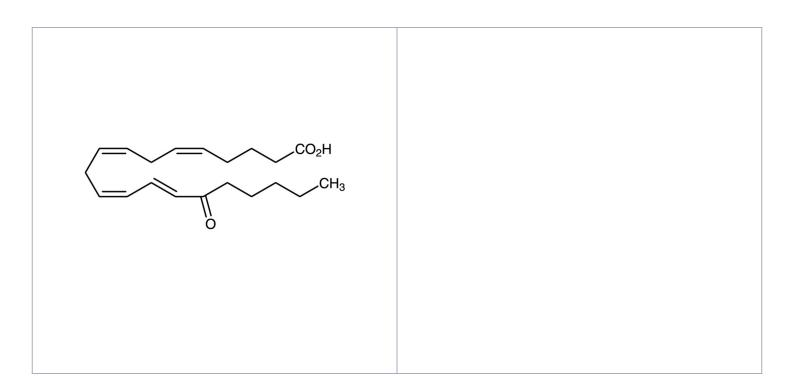
Ordering Information

Order Online »

BML-K015-0050 50µg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Store, as supplied,

> at -80°C for up to 1 year. Product is subject to oxidation. Protect from air. Open ampule and aliquot into suitable tubes or vials. For experiments with cells or tissues, aliquots of ethanol solutions may be added directly to aqueous media. For HPLC, aliquots of ethanol solutions should be evaporated to dryness under a stream of nitrogen and

HPLC mobile phase added to the desired concentration.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Liquid. **Appearance**

CAS 81416-72-0

MW 318.5

Purity ≥95% (HPLC)

Last modified: May 29, 2024



info-